



Automation performance, now for labs of any size.

Compact. Fast-Moving. User-Friendly. Intelligent.

The AssayMate™ workstation is a benchtop-sized, automated liquid handler designed to optimize a wide range of laboratory applications. It integrates technologies such as collision detection, error handling, and real-time run visualization. The software, featuring an intuitive drag-and-drop interface, simplifies protocol building, making automation performance accessible to all.

By streamlining high-volume, repetitive tasks, the AssayMate workstation significantly increases laboratory throughput and enables reproducibility and traceability. This operational efficiency frees up valuable time, allowing scientists to focus on data analysis and exploring new frontiers in scientific discovery.

Some key advantages:

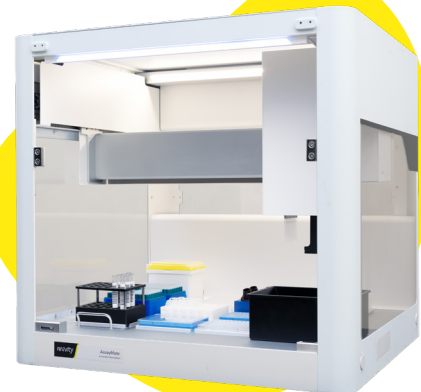
- Partial tip loading capability allows to pipette from tube/vial/bottle
- Reconfigurable recessed deck easily accommodates optional modules
- Wide volume range without needing to switch head

8-Channel	Performance*			
Pipette volume	3-1000 µL			
Pipetting accuracy and precision (Single transfer)	Tip / µL	Dispense	Accuracy	Precision
	50	3 µL	≤6%	≤2%
	50	5 µL	≤2%	≤2%
	200	50 µL	≤2%	≤1%
Feature	Automated pick up, eject and return tip, liquid level detection, liquid level tracking			
Liquid handling	Multi-dispense, multi-aspirate, air gap, aspirate and dispense delay, mix, etc.			

*Performance measured using distilled water and Revvity's pipette tips.

Key features:

- Two configurations:
 - 8 channels + gripper
 - 1+8 pipetting channels
- Air displacement pipetting technology
- Pressure-based liquid level detection (pLLD)
- 12 + 4 (staging slots)
- Suitable for SBS labware
- Compact size
- Budget friendly
- Simple, intuitive software interface
- Optional modules enable many varied applications



| AssayMate™ workstation

Sample prep to accelerate your workflow

Hardware key features

Air displacement pipetting

Two pipetting heads are available: 1 or 8 fixed channels. The 8-channel pipetting head enables partial tip loading, automating medium throughput workflows.

Design efficiency

With 12+4 staging slots and an adaptive force control gripper, the AssayMate system can handle both hard and soft plates with ease.

Operational simplicity

A five-color indicator front light provides real-time system status updates, while its low noise level minimizes distractions in the lab. The deck includes a reconfigurable, recessed section to accommodate optional modules.

Many varied experimental processes can be performed when adding the optional Heater Shaker, Heater Cooler, and Magnetic Block modules.

Software key features

The AssayPREP™ software is built for efficiency and ease of use, with minimal training required.

Intuitive protocol building and editing

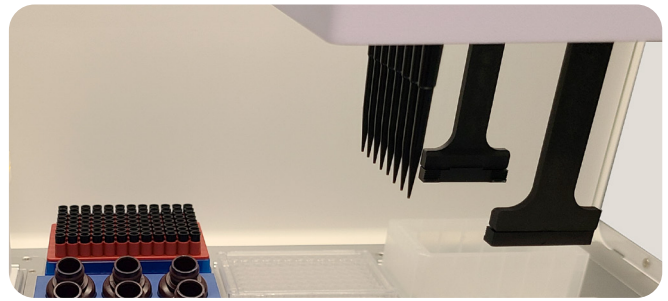
the drag-and-drop interface simplifies the creation of simple to more complex protocols.

Reliable operation

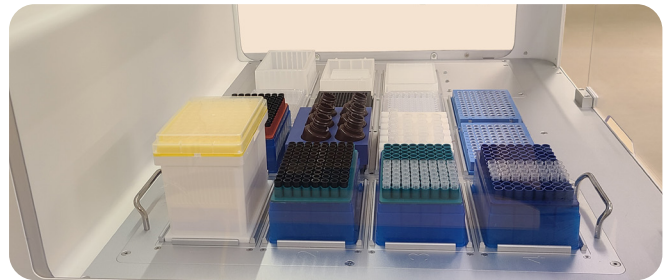
Features like plate drop detection and error handling increase the integrity of your runs while improving safety.

Full protocol visualization

Real-time run visualization provides valuable oversight of your experiment.



| 8-channel pipette and labware gripper



| Modular deck arrangement

Drag-and-drop editing: With drag-and-drop functionality and smart editing suggestions, with no code programming

Easy settings: Support default and custom labware and liquid classes

One-click start: Start experiments immediately and display the experiment process in real-time

Quick recovery: Reliable error handling mechanism* allows experiment resumption

* For most error conditions



Learn more

revvity